

Separation of some nucleotide units

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F254s (P/N 1. 13192, alt P/N 1. 15647)
Eluent	Methanol/water 90/10 (v/v) with addition of 0,2 mol/l lithium chloride
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. ApApA 2. ApA 3. Adensoine (all 0,1%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

